



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human Insulin Kit, SECTOR**Lot Number:** K0082838**Expiration Date:** 30 JUN 2027**Catalog Number:** K151S5-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description		Lot Number	Storage Temperature	Expiration Date
Human Insulin 96-Well Small Spot SECTOR Plate		Z0048201	2-8°C	30 APR 2028
Human Insulin Antibody		D0082539	2-8°C	30 JUN 2029
Metabolic Panel 1 Calibrator Blend		A0080438	2-8°C	31 JUL 2030
Diluent 11		M0110071	≤ -10°C	31 JUL 2027
Diluent 56		M0560009	≤ -10°C	30 JUN 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Metabolic Panel 1 Control 1		Metabolic Panel 1 Control 2		Metabolic Panel 1 Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Insulin	80% - 120%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%	95% - 135%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	Insulin	0.069	80% - 120%	< 20%	19.5	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Insulin	29.3	µIU/mL

Dilution to top of Curve: 1X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 FEB 2026